

Data Sheet

HEXA

Human, recombinant, C-terminal FLAG-His-tags

Catalog #: 71087

Lot#: 130503-H

Conc.: 0.23 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 150 mM imidazole and 20% glycerol

Stability: >6 months at -80 °C

References:

1. Lemieux, M.J., *et al.*, *J Mol Biol.* 2006;**359**:913-929.
2. Peleg, L., *et al.*, *Biochem Mol Med.* 1995 Apr;**54**(2):126-32.

Description:

Human hexosaminidase A (alpha polypeptide), also known as HEXA, (GenBank Accession No. NM_000520), full length (a.a. 1-529(end)) with C-terminal FLAG-His-tags, MW = 62.5 kDa, expressed in a baculovirus infected Sf9 cell expression system.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

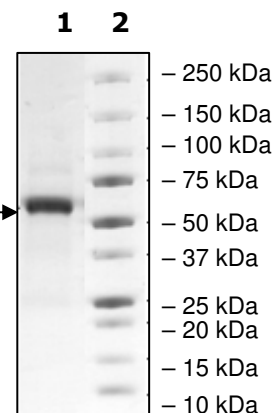
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2.8 µg HEXA

Lane 2:
Protein Marker

MW: 62.5 kDa
Purity: ≥91%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MTSSRLWFSLLLAAAFAGRATALWPWPQNFQTSDQRYVLYPNNFQFQYDVSSAAQPG
CSVLDEAFQRYRDLLFGSGSWPRPYLTGKRHTLEKNLVVSVVTPGCNQLPTLESVEN
YTLTINDDQCLLLSETVWGALRGLETFSQLVWKS AEGTFFINKTEIEDFPRFPHRGLLLD
TSRHYP LSSILDTLDVMAYNKLNVFHWHLVDDPSFPYESFTFPELMRKGSPVTHIYT
AQDVKEVIEYARLRGIRVLA EFDTPGHTLSWGP GIPGLLTPCYSGSEPSGTFGPVNPSL
NNTYEFMSTFFLEVSSVFPDFYLHLGGDEVDFTCWKS NPEIQDFMRKKGFGEDFKQLE
SFYIQTL LDIVSSYGKGYVVWQEVFDNKVKIQPDTIIQVWREDIPVNYMKELELVTKAGFR
ALLSAPWYLN RISYGPDWKDFYIVEPLAFEGTPEQKALVIGGEACMWGEYVDNTNLVP
RLWPRAGAVAERLWSNKLTSDLTFAYERLSHFRCCELLRRGVQAQPLNVGFCEQEFEQ
TDYKDDDDKHHHHHH

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com